

# PRODUCT INFORMATION

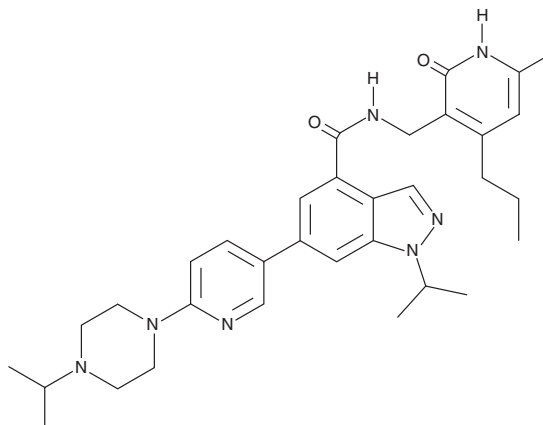


## UNC1999

Item No. 14621

**CAS Registry No.:** 431612-23-5  
**Formal Name:** N-[(1,2-dihydro-6-methyl-2-oxo-4-propyl-3-pyridinyl)methyl]-1-(1-methylethyl)-6-[6-[4-(1-methylethyl)-1-piperazinyl]-3-1H-indazole-4-carboxamide

**MF:** C<sub>33</sub>H<sub>43</sub>N<sub>7</sub>O<sub>2</sub>  
**FW:** 569.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 31, 282, 313 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

UNC1999 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC1999 in the solvent of choice, which should be purged with an inert gas. UNC1999 is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of UNC1999 in these solvents is approximately 0.1 and 1 mg/ml, respectively.

UNC1999 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

The histone H3 lysine 27 (H3K27) methyltransferase EZH2 plays an important role in regulating gene expression, and its aberrant activity is linked to the onset and progression of cancer.<sup>1</sup> UNC1999 is a cell-permeable EZH2 inhibitor (IC<sub>50</sub> = 2 nM) that is 22-fold selective over EZH1 and >1,000-fold selective over other histone methyltransferases.<sup>2</sup> UNC1999 has been shown to inhibit H3K27 methylation in MCF10A cells with an IC<sub>50</sub> value of 124 nM.<sup>2</sup> For more information on UNC1999 please visit the Structural Genomics Consortium (SGC). The negative control, UNC2400, for UNC1999 is also available exclusively through the SGC.

### References

1. Simon, J.A. and Lange, C.A. Roles of the EZH2 histone methyltransferase in cancer epigenetics. *Mutat. Res.* **647(1-2)**, 21-29 (2008).
2. Konze, K.D., Ma, A., Li, F., et al. An orally bioavailable chemical probe of the lysine methyltransferases EZH2 and EZH1. *ACS Chem. Biol.* **8(6)**, 1324-1334 (2013).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/06/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM